



Gene Pulser® Electroprotocols

* We recommend adapting this protocol to use the Gene Pulser electroporation buffer (catalog #165-2676, 165-2677), which increases cell viability and transfection efficiency in mammalian cell lines.

Cell Type	Mammalian, adherent, suspension,	Molecules	DNA: circular, 9.2kB, expression
Species	Monkey, Vero, kidney cells; Human,	Electroporated	vectors
Used	C-4I, cervical carcinoma cells		

Before the Pulse

Cell growth medium	M199 (GIBCO/BRL, Sigma)	Growth phase at harvest	3 x 10 ⁽⁶⁾ cells / ml
Wash solution	M199	Pre-pulse incubation	None

The Pulse

Electroporation Temperature	Room temperature	Instruments Used	Gene Pulser® apparatus & Capacitance Extender
Electroporation Medium*	M199	Cuvette Gap	0.4cm
Cell Density	3 x 10 ⁽⁶⁾ cells / ml	Voltage	0.250 kV
Volume of Cells	0.3 ml / pulse	Field Strength	0.625 kV/cm
DNA Concentration	100 µg DNA / pulse	Capacitor	500 µF
DNA Resuspension Buffer	distilled water	Resistor	(Pulse Controller) Ω none
Volume of DNA	5 µl	Time Constant	Not given

After the Pulse

Outgrowth Medium	M199 - DMEM /F12,5% FBS,5% serum supplement, 1% pen/ strep, 400 µg G418
Outgrowth Temperature	37°C
Length of Incubation	24 hours recovery before selection
Selection Method or Assay Used	G418
Electroporation Efficiency	1.3 / µg DNA
Per Cent Survival	Not given

Relevant Publications and/or Comments

Note: exponential values designated in parentheses.

Name of Submitter
Institution Address

Telephone Number
Fax Number
Date Submitted 8/24/91
Survey Number 155
© Bio-Rad Laboratories, 1993